

## NB100-181 Protocol

# Western Blot protocol for DMC1 Antibody (NB100-181)

DMC1 Antibody (1D12/4): [https://www.novusbio.com/products/dmc1-antibody-1d12-4\\_nb100-181](https://www.novusbio.com/products/dmc1-antibody-1d12-4_nb100-181)

Western Blot

Procedure Guide for NB 100-181 Anti-hDMC-1

Western Blot

1. Proteins are separated on an SDS-PAGE gel.
2. Following the protein transfer, the membrane is blocked with 5% NFDM in TBS-T.
3. The membrane is then washed.
4. NB 100-181 [anti-hDMC] is diluted 1:1,000 in dilution buffer [5% NFDM in TBS-T] and incubated for 3 hours at room temperature, gently shaking.
5. The membrane is then washed.
6. Secondary antibody [HRP conjugated goat anti-mouse] is incubated for 45 minutes at room temperature, gently shaking.
7. Develop membrane.

NOTE: Mouse testis tissue extracts were used as a positive control for this antibody. MEF 161 and HeLa whole cell extracts were not shown to work in Western analysis.